

Product Number:	101419
Lot Number:	501460
Expiration:	20-Feb-19
Platform(s):	HD-1



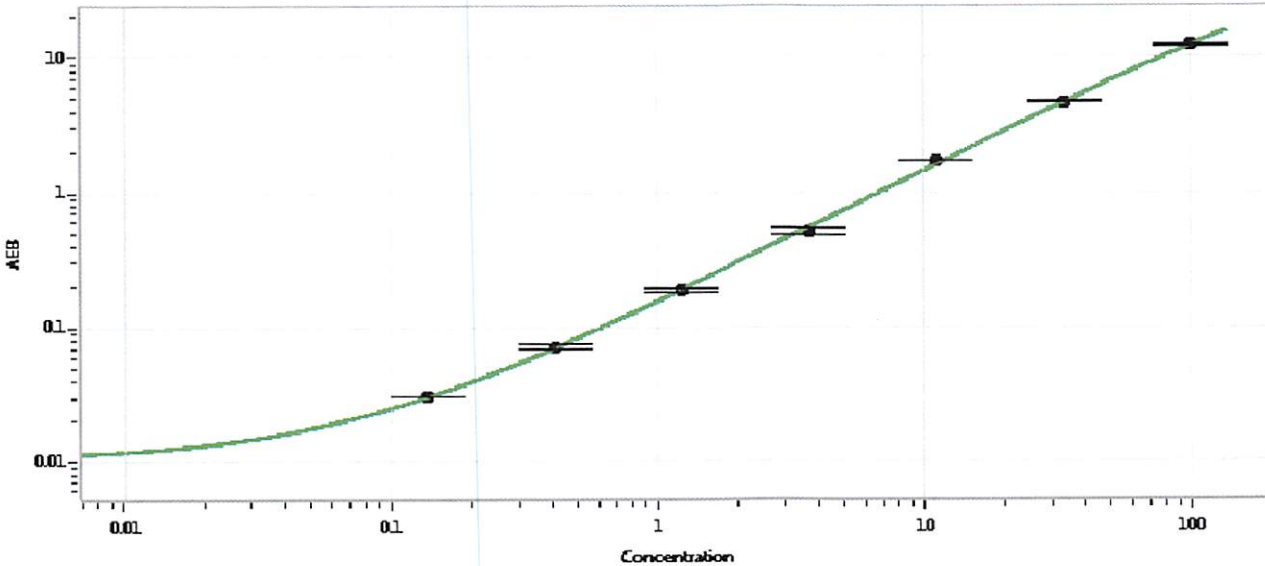
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101899	827602
Detector Stock	101900	828602
Calibrator Concentrate	101420	828601
Calibrator Diluent	101418	823615
Sample Diluent	101416	827407

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
30 OCT 2018